

Dog Liver X Receptor, Alpha (dLXRα, nr1h3) Reporter Assay System

96-well Format Assays Product # D00311

Technical Manual (version 7.1i)

www.indigobiosciences.com

3006 Research Drive, Suite A1, State College, PA 16801, USA

Customer Service: 814-234-1919; FAX 814-272-0152 customerserv@indigobiosciences.com

Technical Service: 814-234-1919 techserv@indigobiosciences.com



Dog LXRa Reporter Assay System 96-well Format Assays

I. Description	
■ The Assay System	3
■ The Assay Chemistry	3
Preparation of Test Compounds	4
• Considerations for Automated Dispensing	4
Assay Scheme	4
Assay Performance	5
II. Product Components & Storage Conditions	6
III. Materials to be Supplied by the User	6
IV. Assay Protocol	
A word about <i>Antagonist</i> -mode assay setup	7
■ DAY 1 Assay Protocol	7
■ DAY 2 Assay Protocol	9
V. Related Products	10
VI. Limited Use Disclosures	11
APPENDIX 1: Example Scheme for Serial Dilution	12

I. Description

The Assay System

This nuclear receptor assay utilizes proprietary non-Dog mammalian cells engineered to provide constitutive, high-level expression of **Dog Liver X Receptor Alpha** (nr1h3), a ligand-dependent transcription factor referred to as $dLXR\alpha$.

The N-terminal DNA binding domains (DBD) of the native LXR α has been substituted with that of the yeast GAL4-DBD. The reporter gene is beetle luciferase functionally linked to the GAL4 upstream activation sequence (UAS). Thus, quantifying changes in luciferase expression in the treated reporter cells provides a specific and sensitive surrogate measure of the changes in LXR α activity. The principal application of this assay is in the screening of test samples to quantify any functional activity, either agonistic or antagonistic, that they may exert against Dog LXR α .

INDIGO's Reporter Cells are prepared using INDIGO's proprietary **CryoMite**TM process. This cryo-preservation method yields exceptional cell viability post-thaw, and provides the convenience of immediately dispensing healthy, division-competent reporter cells into assay plates. There is no need for cumbersome intermediate treatment steps such as spin-and-rinse of cells, viability determinations, cell titer adjustments, or the pre-incubation of reporter cells prior to assay setup.

INDIGO Bioscience's assays are all-inclusive cell-based assay systems. In addition to LXR α Reporter Cells, this kit provides two optimized media for use during cell culture and in diluting the user's test samples, the reference agonist TO901317, Luciferase Detection Reagent, and a cell culture-ready assay plate.

The Assay Chemistry

INDIGO's assays capitalize on the extremely low background, high-sensitivity, and broad linear dynamic range of bio-luminescence reporter gene technology.

Reporter Cells incorporate the cDNA encoding beetle luciferase, a 62 kD protein originating from the North American firefly (*Photinus pyralis*). Luciferase catalyzes the mono-oxidation of D-luciferin in a Mg⁺²-dependent reaction that consumes O₂ and ATP as co-substrates, and yields as products oxyluciferin, AMP, PP_i, CO₂, and photon emission. Luminescence intensity of the reaction is quantified using a luminometer and is reported in terms of Relative Light Units (RLU's).

INDIGO's assays feature a luciferase detection reagent specially formulated to provide stable light emission between 5 and 90+ minutes after initiating the luciferase reaction. Incorporating a 5-minute reaction-rest period ensures that light emission profiles attain maximal stability, thereby allowing assay plates to be processed in batch. By doing so, the signal output from all sample wells, from one plate to the next, may be directly compared within an experimental set.

Preparation of Test Compounds

Small molecule test compounds are typically solvated in DMSO at high concentrations; ideally 1,000x-concentrated stocks relative to the highest desired treatment concentration in the assay. Using high-concentration stocks minimizes DMSO carry-over into the assay plates. Immediately prior to setting up an assay, the master stocks are serially diluted using one of two alternative strategies:

1.) As described in *Step 2* and depicted in Appendix 1 for the reference agonist TO901317, **Compound Screening Medium (CSM)** may be used as the diluent to make serial dilutions of test compounds to achieve the desired final assay concentration series.

Alternatively, if test compound solubility is expected to be problematic,

2.) DMSO may be used to make serial dilutions, thereby generating 1,000x-concentrated stocks for each independent test concentration. Treatment media are then prepared using CSM to make final 1,000-fold dilutions of the prepared DMSO dilution series.

Regardless of the dilution method used, the final concentration of total DMSO carried over into assay wells should not exceed 0.4%. DMSO-induced cytotoxicity can be expected above 0.4%.

NOTE: CSM is formulated to help stabilize hydrophobic test compounds in the aqueous environment of the assay mixture. Nonetheless, high concentrations of extremely hydrophobic test compounds diluted in CSM may lack long-term stability and/or solubility, especially if further stored at low temperatures. Hence, it is recommended that test compound dilutions are prepared in CSM immediately prior to assay setup and are considered to be 'single-use' reagents.

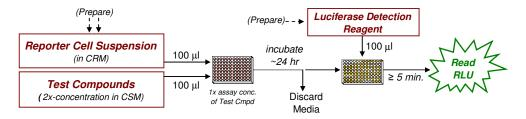
Considerations for Automated Dispensing

When processing a small number of assay plates, first carefully considered the dead volume requirement of your dispensing instrument before committing assay reagents to its setup. In essence, "dead volume" is the volume of reagent that is dedicated to the instrument; it will *not* be available for final dispensing into assay wells. The following Table provides information on reagent volume requirements, and available excesses.

Stock Reagent & Volume provided	Volume to be Dispensed (96-well plate)	Excess rgt. volume available for instrument dead volume
Reporter Cell Suspension 12 ml (prepared from kit components)	100 μl / well 9.6 ml / plate	~ 2.4 ml
LDR 12 ml (prepared from kit components)	100 μl / well 9.6 ml / plate	~ 2.4 ml

Assay Scheme

Figure 1. Assay workflow. *In brief*, Reporter Cells are dispensed into wells of the assay plate and then immediately dosed with the user's test compounds. Following 22 -24 hours incubation, treatment media are discarded and prepared Luciferase Detection Reagent (LDR) is added. Light emission from each assay well is quantified using a plate-reading luminometer.



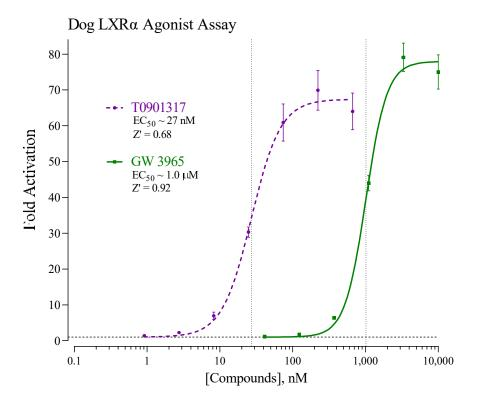


Figure 2. Dog LXRα agonist dose-response assays.

Dog LXR α Reporter Cells were treated with the reference agonists T0901317 (provided) and GW3965 (Tocris). Luminescence was quantified and values of Average RLU and corresponding standard deviation and percent coefficient of variation, and Fold-Activation were calculated for each treatment group (n = 4). Z' values were calculated as described by Zhang, *et al.* (1999)\(^1\). GraphPad Prism software was used to perform the least-squares method of non-linear regression to curve-fit data and determine EC50 values.

$$Z' = 1 - [3*(SD^{Ref} + SD^{Bkg.}) / (RLU^{Ref.} - RLU^{Bkg.})]$$

¹ Zhang JH, Chung TD, Oldenburg KR. (1999) A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. J Biomol Screen.:**4**(2), 67-73.

II. Product Components & Storage Conditions

This Dog LXR α assay kit contains materials to perform assays in a single 96-well assay plate.

Reporter cells are temperature sensitive! To ensure maximal viability the tube of cells must be maintained at -80° C until immediately prior to the rapid-thaw procedure described in *Step 3* of this protocol.

Assay kits are shipped on dry ice. Upon receipt of the kit transfer it to -80°C storage. If you wish to first inspect and inventory the individual kit components be sure to first transfer and submerge the tube of reporter cells in dry ice.

The aliquot of Reporter Cells is provided as a single-use reagent. Once thawed, reporter cells can NOT be refrozen, nor can they be maintained in extended culture with any hope of retaining downstream assay performance. Therefore, extra volumes of these reagents should be discarded after assay setup.

The date of product expiration is printed on the Product Qualification Insert (PQI) enclosed with each kit.

Kit Components	<u>Amount</u>	Storage Temp.
- Dog LXRα Reporter Cells	1 x 2.0 mL	-80°C
• Cell Recovery Medium (CRM)	1 x 10.5 mL	-20°C
• Compound Screening Medium (CSM)	1 x 35 mL	-20°C
• TO901317, 600 μM (in DMSO) (reference agonist for LXR's)	1 x 30 μL	-20°C
Detection Substrate	1 x 6.0 mL	-80°C
• Detection Buffer	1 x 6.0 mL	-20°C
 96-well assay plate (white, sterile, cell-culture ready) 	1	ambient

III. Materials to be Supplied by the User

The following materials must be provided by the user, and should be made ready prior to initiating the assay procedure:

DAY 1

- dry ice bucket (Step 3)
- cell culture-rated laminar flow hood.
- 37°C, humidified 5% CO₂ incubator for mammalian cell culture.
- 37°C water bath.
- 70% alcohol wipes
- 8-channel electronic, repeat-dispensing pipettes & sterile tips
- disposable media basins, sterile.
- sterile multi-channel media basins (such as the Heathrow Scientific "Dual-Function Solution Basin"), *or* sterilized 96 deep-well blocks (*e.g.*, Axygen Scientific, #P-2ML-SQ-C-S), *or* appropriate similar vessel for generating dilution series of reference and test compound(s).
- Optional: antagonist reference compound.
- Optional: clear 96-well assay plate, sterile, cell culture treated, for viewing cells on Day 2.

DAY 2 plate-reading luminometer.

IV. Assay Protocol

Review the entire Assay Protocol before starting. Completing the assay requires an overnight incubation. *Steps 1-8* are performed on *Day 1*, requiring less than 2 hours to complete. *Steps 9-14* are performed on *Day 2* and require less than 1 hour to complete.

A word about Antagonist-mode assay setup

Receptor inhibition assays expose the Reporter Cells to a constant, sub-maximal concentration (typically between $EC_{50} - EC_{85}$) of a known agonist AND the test compound(s) to be evaluated for antagonist activity. This LXR α assay kit includes a 600 μ M stock of **TO901317**, a potent agonist of the LXR's that may be used to setup antagonist-mode assays. 160 nM TO901317 typically approximates EC_{80} in this assay. Hence, it presents a suitable assay concentration of agonist to be used when screening test compounds for inhibitory activity.

Adding a 2x-EC₈₀ concentration (~320 nM) of the challenge agonist TO901317 to the bulk suspension of Reporter Cells (*i.e.*, prior to dispensing into assay wells) is the most efficient and precise method of setting up antagonist assays, and it is the method presented in *Step 5b* of the following protocol. Note that, in *Step 6*, 100 µl of treatment media is combined with 100 µl of pre-dispensed [Reporter Cells + agonist]. Consequently, one must prepare the bulk suspension of Reporter Cells to contain a 2x-EC₈₀ concentration of TO901317. **APPENDIX 1** provides a dilution scheme that may be used as a guide when preparing cell suspension supplemented with a 2x-concentration of TO901317.

DAY 1 Assay Protocol: All steps must be performed using aseptic technique.

- **1.**) Remove **Cell Recovery Medium (CRM)** and **Compound Screening Medium (CSM)** from freezer storage and thaw in a 37°C water bath.
- **2.) Prepare dilutions of treatment compounds** (first see *Note 5.4*): Prepare Test Compound treatment media for *Agonist-* or *Antagonist-mode* screens.

The final concentration of total DMSO, or any organic solvent, carried over into assay reactions should not exceed 0.4%.

Note that, in Step 6, 100 μ l of the prepared treatment media is added into assay wells that have been pre-dispensed with 100 μ l of Reporter Cells. Hence, to achieve the desired *final* assay concentrations one must prepare treatment media with a 2x-concentration of the test and reference material(s). Use **CSM** to prepare the appropriate dilution series. Manage dilution volumes carefully. This assay kit provides 35 ml of CSM.

Preparing the positive control: This LXRα assay kit includes a 600 μ M stock solution of **TO901317**, a potent reference agonist of LXRα. The following 7-point treatment series, with concentrations presented in 3-fold decrements, provides a suitable dose-response: 600, 200, 66.7, 22.2, 7.41, 2.47, and 0.823 nM. Always include 'no treatment' (or 'Vehicle') control wells. **APPENDIX 1** provides an example for generating such a dilution series.

3.) Rapid Thaw of the Reporter Cells: *First*, retrieve the tube of **CRM** from the 37°C water bath and sanitize the outside surface with a 70% ethanol swab.

Second, retrieve **Reporter Cells** from -80°C storage and *immerse in dry ice* to transport the tube to a laminar flow hood. When ready, transfer the tube of cells from dry into a rack and, without delay, perform a *rapid thaw* of the frozen cells by transferring a 10 ml volume of 37°C CRM into the tube of frozen cells. Recap the tube of cells and immediately place it in a 37°C water bath for 5 - 10 minutes. The resulting volume of cell suspension will be 12 ml.

4.) Retrieve the tube of Reporter Cell Suspension from the water bath. Sanitize the outside surface of the tube with a 70% alcohol swab.

5.) *a. Agonist*-mode assays. Gently invert the tube of Reporter Cells three times to gain a homogenous cell suspension. Without delay, dispense $\underline{100 \, \mu l}$ of cell suspension into each well of the assay plate.

~ or ~

- **b.** Antagonist-mode assays. Supplement the bulk suspension of Reporter Cells with the desired $\underline{2x}$ -concentration of TO901317 (refer to "A word about antagonist-mode assay setup", pg. 7). Gently invert the tube of Reporter Cells three times to gain a homogenous cell suspension. Dispense 100 μ l of cell suspension into each well of the assay plate.
 - NOTE 5.1: If INDIGO's Live Cell Multiplex Assay is to be incorporated, a minimum of 3 'cell blank' wells (meaning cell-free, but containing 'Compound Screening Media') must be included in the assay plate to allow quantification of plate-specific fluorescence background (refer to the LCMA Technical Manual).
 - *NOTE 5.2:* Take special care to prevent cells from settling during the dispensing period. Allowing cells to settle during the transfer process, and/or lack of precision in dispensing uniform volumes across the assay plate *will* cause well-to-well variation (= increased Standard Deviation) in the assay.
 - NOTE 5.3: Users sometimes wish to examine the reporter cells using a microscope. If so, the extra volume of cell suspension provided with each kit may be dispensed (100 μ l/well) into a clear 96-well cell culture treated assay plate, followed by 100 μ l/well of CSM. Incubated overnight in identical manner to those reporter cells contained in the white assay plate.
 - *NOTE 5.4:* For logistical reasons, some users find it more convenient to first plate the reporter cells and then prepare their test compound dilutions. That strategy works equally well. Once plated, cells may be placed in an incubator for up to 3 hours before proceeding to *Step 6*.
 - *NOTE 5.5:* If well-to-well variation due to 'edge-effects' is a concern this problem *may* be mitigated by dispensing sterile liquid into the *inter-well* spaces of the assay plate. Simply remove 1 tip from the 8-chanel dispenser and dispense 100 μ l of sterile water into each of the seven inter-well spaces per column of wells.
- **6.)** Dispense 100 µl of 2x-concentration treatment media into appropriate assay wells.
- **7.)** Transfer the assay plate into a cell culture incubator (37°C, humidified 5% CO₂) for 22 24 hours.
 - *NOTE:* Ensure a high-humidity ($\geq 70\%$) environment within the cell culture incubator. This is critical to prevent the onset of deleterious "edge-effects" in the assay plate.
- **8.)** For greater convenience on *Day 2*, retrieve **Detection Substrate** *and* **Detection Buffer** from freezer storage and place them in a dark refrigerator (4°C) to thaw overnight.

- **DAY 2 Assay Protocol:** Subsequent manipulations do *not* require special regard for aseptic technique and may be performed on a bench top.
- **9.**) 30 minutes before intending to quantify receptor activity, remove **Detection Substrate** and **Detection Buffer** from the refrigerator and place them in a low-light area so that they may equilibrate to room temperature.
 - *NOTE:* Do NOT actively warm Detection Substrate above room temperature. If these solutions were not allowed to thaw overnight at 4°C, a room temperature water bath may be used to expedite thawing.
- **10.**) Set the plate-reader to "luminescence" mode. Program the instrument to perform a single <u>5 second</u> "plate shake" prior to reading the first assay well. Set read-time to 0.5 second (500 mSec) per well, *or less*.
- **11.**) Immediately before proceeding to *Step 12*, prepare **Luciferase Detection Reagent** (**LDR**). Combine 'Detection Buffer' and 'Detection Substrate' by pouring-over their entire volumes into a media basin; rock the basin gently to mix the reagent. The resulting volume of LDR is 12 ml.
- **12.**) Following 22 24 hours of incubation discard all media contents. The preferred method is to use a 'wrist flick' to eject the treatment media into an appropriate waste container. *Gently* tap the inverted plate onto a clean absorbent paper towel to remove residual droplets. Cells will remain tightly adhered to well bottoms.
- 13.) Add $\underline{100 \,\mu l}$ of **LDR** to each well of the assay plate. Allow the assay plate to rest at room temperature for at least $\underline{5 \, \text{minutes}}$. Do not shake the assay plate during this period.
- 14.) Quantify luminescence.

V. Related Products

Product No.	Product Descriptions		
Human LXRα Assay Kits			
IB00311-32	Human LXRα Reporter Assay System 3x 32 assays in 8-well strips, 96-well format		
IB00311	Human LXRα Reporter Assay System 1x 96-well format assay		
IB00312	Human LXRα Reporter Assay System 1x 384-well format assays		
Bulk volumes of Assay Reagents may be custom manufactured to accommodate any scale of HTS. Please Inquire.			
Dog LXRα Assay Kits			
D00311	Dog LXRα Reporter Assay System 1x 96-well format assay		
	Mouse LXRα Assay Kits		
M00311-32	Mouse LXRα Reporter Assay System 3x 32 assays in 8-well strips, 96-well format		
M00311	Mouse LXRα Reporter Assay System 1x 96-well format assay		
Rat LXRα Assay Kits			
R00311-32	Rat LXRα Reporter Assay System 3x 32 assays in 8-well strips, 96-well format		
R00311	Rat LXRα Reporter Assay System 1x 96-well format assay		

LIVE Cell Multiplex (LCM) Assay		
LCM-01	Reagent volumes sufficient to perform 96 Live Cell Assays in 1x96-well, or 2x48-well, or 3x32-well assay plate formats	
LCM-05	Reagent in 5x bulk volume to perform 480 Live Cell Assays performed in 5 x 96-well assay plates	
LCM-10	Reagent in 10x bulk volume to perform 960 Live Cell Assays performed in 10 x 96-well assay plates	
INDIGIo Luciferase Detection Reagent		
LDR-10, -25, -50, -500	INDIGIo Luciferase Detection Reagents in 10 mL, 25 mL, 50 mL, and 500 mL volumes	

Please refer to INDIGO Biosciences' website for updated product offerings.

www.indigobiosciences.com

VI. Limited Use Disclosures

Products commercialized by INDIGO Biosciences, Inc. are for RESEARCH PURPOSES ONLY – not for therapeutic, diagnostic, or contact use in humans or animals.

"CryoMite" is a Trademark TM of INDIGO Biosciences, Inc. (State College, PA, USA).

Product prices, availability, specifications, claims and technical protocols are subject to change without prior notice. The printed Technical Manual provided in the kit box will always be the most recently updated version.

Copyright © INDIGO Biosciences, Inc. (State College, PA, USA). All rights reserved.

APPENDIX 1

Example scheme for the serial dilution of TO901317 reference agonist and the setup of an LXR α dose-response assay.

